

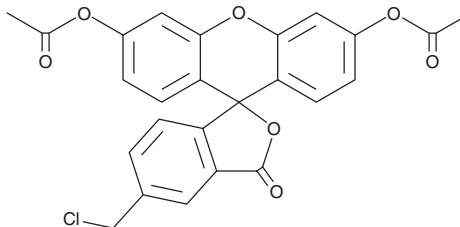
PRODUCT INFORMATION



Green CMFDA

Item No. 19583

CAS Registry No.: 136832-63-8
Formal Name: 3',6'-bis(acetyloxy)-5-(chloromethyl)-spiro[isobenzofuran-1(3H),9'-[9H]xanthen]-3-one
Synonyms: 5-Chloromethylfluorescein diacetate
MF: C₂₅H₁₇ClO₇
FW: 464.9
Purity: ≥95%
UV/Vis.: λ_{max}: 220, 291 nm
Ex./Em. Max: 492/517 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Green CMFDA is supplied as a crystalline solid. A stock solution may be made by dissolving the green CMFDA in the solvent of choice. Green CMFDA is miscible in the organic solvent DMSO, which should be purged with an inert gas.

Description

Green CMFDA is a cell-permeable fluorescent probe that freely enters living cells, where it is transformed into a cell-impermeable, soluble, fluorescent compound. The resulting molecule does not affect cell viability or proliferation and stably fluoresces for at least 72 hours, allowing for its use in cell tracking.¹ It is also transferred to daughter cells. Green CMFDA is brightly fluorescent with standard filters (ex/em max = 492/517 nm) and is often used with additional fluorescent dyes and proteins.^{2,3}

References

1. Beem, E., and Segal, M. S. Evaluation of stability and sensitivity of cell fluorescent labels when used for cell migration. *J. Fluoresc.* **23(5)**, 975-987 (2013).
2. Dagley, M. J., Saunders, E. C., Simpson, K. J., et al. High-content assay for measuring intracellular growth of *Leishmania* in human macrophages. *Assay Drug Dev. Technol.* **13(7)**, 389-401 (2015).
3. Elliott, J. T., Tona, A., and Plant, A. L. Comparison of reagents for shape analysis of fixed cells by automated fluorescence microscopy. *Cytometry* **52(2)**, 90-100 (2003).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/13/2017

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM